



EDA-GTPyS-ATTO-655

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO 655, Triethylammonium salt

Cat. No.	Amount
NU-1610-655	50 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤ 2 ml).

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₄₀H₅₃N₁₀O₁₉P₃S₂ (free acid)

Molecular Weight: 1134.96 g/mol (free acid)

Exact Mass: 1134.21 g/mol (free acid)

Purity: ≥ 90 % (HPLC)

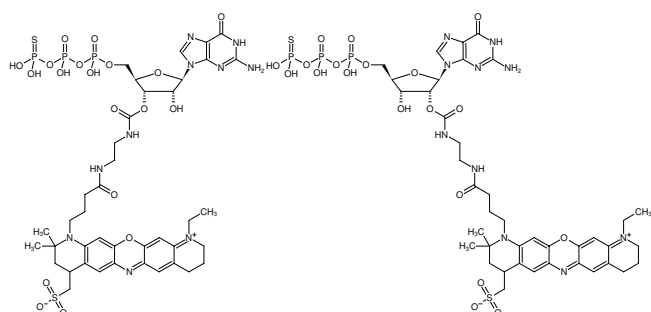
Form: solution in water

Color: blue

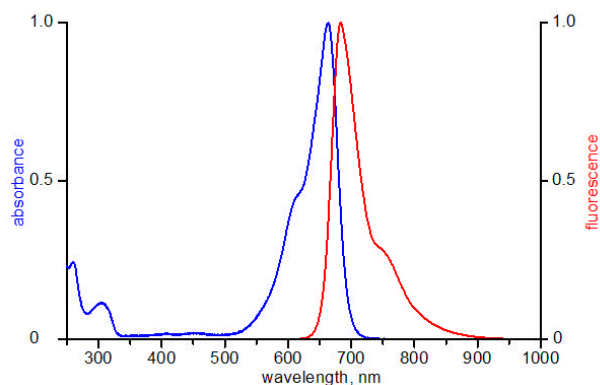
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 663 nm, λ_{em} 680 nm, ε 125.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-GTPyS-ATTO-655



excitation and emission spectrum of ATTO 655